

Glutathione S- Transferase monoclonal antibody (8C1-G4)

The GST superfamily of enzymes play an important role in detoxification of xenobiotics by catalyzing the conjugation of many hydrophobic and electrophilic compounds with reduced glutathione. There are four main GST classes; alpha, mu, pi, and tau. GST-P is a member of the pi class of GSTs, and its expression is specifically increased in the early stages of rat hepatocarcinogenesis. Due to its high affinity for glutathione, GST is frequently utilized as an approximately 26 kDa fusion tag for the expression and purification of recombinant proteins.

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name GST

Application ELISA, IP, WB

Clone 8C1-G4

Formulation Liquid. In PBS containing 0.01% sodium azide.

Host Mouse

Immunogen GST fusion protein.

Isotype IgG2b

Purity Detail Protein G affinity purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Species independent

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)